

Product Information Sheet

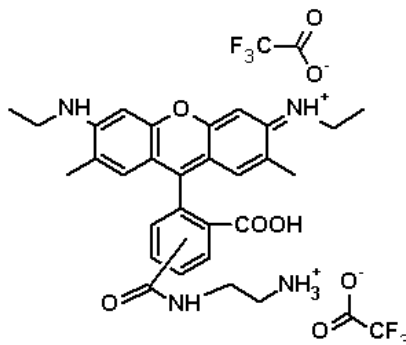
Ordering Information

Product Number: 344
Product Name: 5(6)-Caroxyrhodamine 6G ethylenediamine
Unit Size: 1 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 728.63

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 524 nm
Emission Wavelength: 551 nm

Application Notes

5(6)-Caroxyrhodamine 6G ethylenediamine is a building block for developing rhodamine bioconjugates. In general, rhodamine 6G conjugates have higher fluorescence quantum yields than tetramethylrhodamine conjugates, as well as good photostability. The excitation and emission wavelengths of rhodamine 6G fall between those of fluorescein and tetramethylrhodamine derivatives, making 5(6)-CR6G another color choice for the multicolor fluorescence imaging applications. 5(6)-CR6G is the mixture of two carboxy rhodamine 6G isomers.